

STA13 Antibody [#C11950]

Catalog Number:	C11950
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y3M8
Other Names:	STA13, START domain containing 13, START domain- containing protein 13, StAR-related lipid transfer protein 13, deleted in liver cancer 2, rho GTPase activating protein on chromosome 13q12
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human STA13.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	STA13 antibody detects endogenous levels of total STA13 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:10000
References:	Ching Y.-P., J. Biol. Chem. 278:10824-10830(2003). Nagaraja G.M., Biochem. Biophys. Res. Commun. 313:654-665(2004). Leung T.H.-Y., Proc. Natl. Acad. Sci. U.S.A. 102:15207-15212(2005).